

Product Information

5-Tetramethylrhodamine-dUTP, 1 mM in pH 7.5 Tris-HCl buffer

Catalog Number 40001

Unit Size 25 μ L

Molecular Weight 990.5

Color and Form Red solution

Absorption/Emission 553/577 nm

Storage and Handling

Store Tetramethylrhodamine-dUTP at -20°C or -70°C , protected from light. When stored as recommended, product is stable for at least 6 months from date of receipt.

Product Application

Fluorescent conjugates of dUTP can be used for detection of apoptotic cells by direct fluorescence TUNEL labeling of DNA strand breaks in cells¹, and can be used in place of dTTP in standard DNA labeling and synthesis protocols to generate fluorescent dsDNA and oligonucleotide probes.

Note: The product is not compatible to PCR application.

References

1. Gold et al. (1994). *Lab Invest.* 71 (2):219-25.
2. Slupphaug et al. (1993). *Anal Biochem.* 211 (1):164-9.
3. Hogrefe et al. (2002). *PNAS* 99 (2): 596-601.

Please visit our website at www.biotium.com to view our full selection of CF™ dye bioconjugates, including antibodies, antibody labeling kits, phalloidin, annexin V and a-bungarotoxin, as well as classic fluorescent nucleic acid dyes and antifade fluorescence mounting media.